

**Purified Mouse Anti-human/pig/cow
Cytokeratin 8 Antibody *C-51, monoclonal,
Cross Adsorbed***

Catalog number: V103510
Unit size: 0.1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, pig, cow, sheep
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Cytokeratin 8
Clone	C-51

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse, rabbit, dog, rat, xenopus, chicken serum and then conjugated with under optimal conditions
Application	IHC(P), IP, WB, ICC

Applications

Keratin, type II cytoskeletal 8 is a 54 kDa protein that can be located in the cytoplasm, dystrophin-associated glycoprotein complex and costamere of cells. It is sometimes referred to as keratin-8, type-II keratin Kb8 and cytokeratin-8. Sequencing of keratin, type II cytoskeletal 8 has shown evidence of it contains a primary structural unit, the IF rod domain. Keratin, type II cytoskeletal 8 binds to protein-containing complex and scaffold protein. It is the subject of extensive study due to the fact that it is a member of the extrinsic apoptotic signaling pathway and tumor necrosis factor-mediated signaling pathway. It is an integral part of organismal processes, for example, keratinization, cornification and response to other organism. Mutations and abnormalities in Keratin, type II cytoskeletal 8 have been thought to be involved with a number of diseases, in particular, cirrhosis.